

HSV-2 gG (residues 525-578)

Herpes Simplex Virus-2 glycoprotein G
recombinant, *E. coli*

Cat. No.	Amount
PR-1111-1	1 mg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 25 mM Tris-HCl pH 8.0, 5 mM glutathione, 1.5 M urea and 50% glycerol.

Protein synonyms/aliases

Glycoprotein G.

Application

Recombinant rHSV-2 gG Antigen may be used in ELISA and Western blots, excellent for detection of HSV with minimal specificity problems.

Specificity

Immunoreactive with sera of HSV-infected individuals.

MW

about 35 kDa

Tag

GST

Purity

>95% by SDS-PAGE

Description

Recombinant HSV-2-gG contains the immunodominant regions of glycoprotein G, amino acids: 525-578. HSV-2-gG is purified by proprietary chromatographic techniques.

Background

Glycoprotein G in HSV-2 is the only HSV envelope protein to be cleaved posttranslationally during processing. In the virus-infected cell, gG-2 is cleaved into a secreted amino-terminal portion (sgG-2) and a carboxy-terminal portion. The latter protein is further O-glycosylated, generating the cell membrane-associated mature gG-2 (mgG-2). The mgG-2 protein has widely been used as a prototype antigen for detection of type-specific antibodies against HSV-2.

HSV-2 gG (residues 525-578)

Herpes Simplex Virus-2 glycoprotein G
recombinant, *E. coli*

Selected References:

Morrow *et al.* (2004) Development and use of a type-specific antibody avidity test based on herpes simplex virus type 2 glycoprotein G. *Sex. Transm. Dis.* **31**:508.

Bugli *et al.* (2004) Human monoclonal antibody fragment specific for glycoprotein G in herpes simplex virus type 2 with applications for serotype-specific diagnosis. *J. Clin. Microbiol.* **42**:1250.

Morrow *et al.* (2003) Performance of the focus and Kalon enzyme-linked immunosorbent assays for antibodies to herpes simplex virus type 2 glycoprotein G in culture-documented cases of genital herpes. *J. Clin. Microbiol.* **41**:5212.

Gorander *et al.* (2003) Secreted portion of glycoprotein g of herpes simplex virus type 2 is a novel antigen for type-discriminating serology. *J. Clin. Microbiol.* **41**:3681.

Ikoma *et al.* (2002) Use of a fragment of glycoprotein G-2 produced in the baculovirus expression system for detecting herpes simplex virus type 2-specific antibodies. *J. Clin. Microbiol.* **40**:2526.

Liljeqvist *et al.* (2002) Monoclonal antibodies and human sera directed to the secreted glycoprotein G of herpes simplex virus type 2 recognize type-specific antigenic determinants. *J. Gen. Virol.* **83**:157.